

Product Information

Description

Thalidomide-O-PEG4-NHS ester is a synthesized compound that incorporates the Thalidomide based cereblon ligand and 4-unit PEG linker used in PROTAC technology it is reactive with amine moiety.

Molecular Formula:

$C_{28}H_{33}N_3O_{13}$

Supplied As:

Solid

Solubility:

DCM

Stability:

At least 6 months at -20°C. Avoid freeze/thaw cycles.

Storage:

-20°C

MW:

619.6 Da

Cas No.

2411681-88-2

Purity:

98%

Chemical Structure

